

Guanosine Nucleotide Library

Cat. No.
LIB-102

Purity
> 95% (HPLC)

Storage
-20 °C

Shelf life
12 months

Description

The **Guanosine Nucleotide Library** comprises 80 Guanosine nucleotides, delivered in 1.4 ml size polypropylene tubes packed in 96 well racks, each containing 200 µl of a 1 mM aqueous solution at pH 7.5. Columns 11 and 12 are left empty for control purposes.

Composition

Number	Nucleotide	Cat. No.	Number	Nucleotide	Cat. No.
A 1	GMP, sodium salt	NU-1028	C 1	6-Thio-GDP, sodium salt	NU-1120
A 2	GDP, sodium salt	NU-1172	C 2	6-Thio-GTP, sodium salt	NU-1106
A 3	GTP, sodium salt	NU-1012	C 3	6-Methylthio-GMP, sodium salt	NU-1128
A 4	dGMP, sodium salt	NU-1029	C 4	6-Methylthio-GDP, sodium salt	NU-1129
A 5	dGDP, sodium salt	NU-1030	C 5	6-Methylthio-GTP, sodium salt	NU-1130
A 6	dGTP, sodium salt	NU-1003	C 6	8-Oxo-GTP, sodium salt	NU-1116
A 7	cGMP, sodium salt	NU-1501	C 7	8-Oxo-dGMP, sodium salt	NU-1153
A 8	ddGTP, sodium salt	NU-1017	C 8	8-Oxo-dGDP, sodium salt	NU-1158
A 9	7-Deaza-dGMP, sodium salt	NU-1177	C 9	8-Oxo-dGTP, sodium salt	NU-1117
A 10	7-Deaza-dGDP, sodium salt	NU-1178	C 10	GMP α S, sodium salt	NU-1160
A 11	empty		C 11	empty	
A 12	empty		C 12	empty	
B 1	7-Deaza-dGTP, sodium salt	NU-1179	D 1	GTP α S, sodium salt	NU-409
B 2	2',5'-pGp, sodium salt	NU-1108	D 2	dGDP α S, sodium salt	NU-1164
B 3	3',5'-pGp, sodium salt	NU-1107	D 3	dGTP α S, sodium salt	NU-424
B 4	3'-dGMP, sodium salt	NU-1143	D 4	GpCp, sodium salt	NU-414
B 5	3'-dGDP, sodium salt	NU-1144	D 5	GpCp ₂ , sodium salt	NU-405
B 6	3'-dGTP, sodium salt	NU-1145	D 6	dGpCp ₂ , sodium salt	NU-431
B 7	m ⁷ GMP, sodium salt	NU-1135	D 7	GppCp, sodium salt	NU-402
B 8	m ⁷ GDP, sodium salt	NU-1134	D 8	GppNp, sodium salt	NU-401
B 9	m ⁷ GTP, sodium salt	NU-1122	D 9	GppNH ₂ , sodium salt	NU-1136
B 10	6-Thio-GMP, sodium salt	NU-1121	D 10	GDP β S, sodium salt	NU-427
B 11	empty		D 11	empty	
B 12	empty		D 12	empty	

Guanosine Nucleotide Library

Cat. No.

LIB-102

Purity

> 95% (HPLC)

Storage

-20 °C

Shelf life

12 months

Number	Nucleotide	Cat. No.	Number	Nucleotide	Cat. No.
E 1	GTP _γ S, lithium salt	NU-412	G 1	8-[(6-Amino)hexyl]-amino-cGMP, sodium salt	NU-832
E 2	AP ₄ G, triethylammonium salt	NU-503	G 2	8-[(6-Amino)hexyl]-amino-GMP, sodium salt	NU-829
E 3	AP ₃ G, triethylammonium salt	NU-504	G 3	8-[(6-Amino)hexyl]-amino-GTP, sodium salt	NU-830
E 4	GP ₂ G, triethylammonium salt	NU-531	G 4	EDA-6-Thio-GTP, sodium salt	NU-1613
E 5	GP ₄ G, triethylammonium salt	NU-533	G 5	EDA-GDP, sodium salt	NU-840
E 6	m ⁷ GP ₃ C, triethylammonium salt	NU-512	G 6	EDA-GTP, sodium salt	NU-820
E 7	m ⁷ GP ₅ G, triethylammonium salt	NU-532	G 7	EDA-m ⁷ GDP, sodium salt	NU-827
E 8	AP ₄ (8)G triethylammonium salt	NU-510	G 8	EDA-m ⁷ GTP, sodium salt	NU-824
E 9	AP ₃ (8)G triethylammonium salt	NU-511	G 9	2'-OMe-GTP, sodium salt	NU-1127
E 10	8I-GDP, sodium salt	NU-123	G 10	mant-GDP, triethylammonium salt	NU-204
E 11	empty		G 11	empty	
E 12	empty		G 12	empty	
F 1	8I-GTP, sodium salt	NU-125	H 1	mant-GTP, triethylammonium salt	NU-206
F 2	8Br-GMP, sodium salt	NU-144	H 2	mant-dGTP, triethylammonium salt	NU-212
F 3	8Br-GDP, sodium salt	NU-124	H 3	mant-GTP _γ S, triethylammonium salt	NU-209
F 4	8Br-GTP, sodium salt	NU-118	H 4	2'-Mant-3'-dGTP, triethylammonium salt	NU-231
F 5	6Cl-GTP, sodium salt	NU-143	H 5	ant-GTP, triethylammonium salt	NU-230
F 6	7-Deaza-7I-GTP, sodium salt	NU-1182	H 6	ant-m ⁷ GMP, triethylammonium salt	NU-229
F 7	γ-[(6-Amino)hexyl]-imido]-GTP, sodium salt	NU-816	H 7	ant-m ⁷ GDP, triethylammonium salt	NU-228
F 8	γ-[(6-Amino)hexyl]-imido]-dGTP, sodium salt	NU-849	H 8	ant-m ⁷ GTP, triethylammonium salt	NU-227
F 9	γ-[(8-Amino)octyl]-imido]-GTP, sodium salt	NU-817	H 9	TNP-GDP, triethylammonium salt	NU-217
F 10	α-[(6-Amino)hexyl]-imido]-dGMP, sodium salt	NU-845	H 10	TNP-GTP, triethylammonium salt	NU-220
F 11	empty		H 11	empty	
F 12	empty		H 12	empty	